



# Liquid Assayed Multiquel<sup>®</sup> Premium

Comprehensive Serum Chemistry Control



## Liquid Assayed Multiquel® Premium

Liquid Assayed Multiquel® Premium is a liquid, human serum based comprehensive chemistry control that includes general chemistry, Immunology and Therapeutic Drug Monitoring (TDM) analytes. This assayed trilevel control contains 96 claimed analytes:

- Includes CRP
- Offers 7 additional Immunology analytes, including Cystatin C
- Contains 8 additional TDM analytes, including Vancomycin
- Offers Bile Acids and GLDH<sup>1</sup>
- 3 year shelf life at -20°C to -70°C
- Thawed and unopened stability for 30 days at 2-8°C for most analytes
- 14 day open-vial stability at 2-8°C for most analytes
- Each level sold separately



<sup>1</sup>GLDH is available outside the United States only.

### Analytes

α-1 Acid Glycoprotein*	Bilirubin (Neonatal)	Cystatin C*	Lactate (Lactic Acid)	Salicylate
α-1-Antitrypsin	Bilirubin (Total)	Digoxin	Lambda Light Chain*	Sodium
β-2-Microglobulin*	Caffeine*	Disopyramide*	LAP Arylamidase	T3 Uptake/T-Uptake
Acetaminophen	Calcium (Ionized)	Ethanol	Lidocaine*	T3 (Total)
Acid Phosphatase	Calcium (Total)	Ferritin	Lipase	T4 (Free)
Albumin	Carbamazepine	Gentamicin	Lithium	T4 (Total)
Alpha-Hydroxybutyrate	Carbon Dioxide (CO <sub>2</sub> )	GGT	Magnesium	Theophylline
Dehydrogenase (αHBDH)	Ceruloplasmin	GLDH*	Methotrexate*	Tobramycin
Alkaline Phosphatase	Chloride	Globulin	NAPA*	Transferrin
ALT/SGPT	Cholesterol (HDL)	Glucose	Osmolality	Triglycerides
Amikacin	Cholesterol (LDL)	Haptoglobin	Phenobarbital	TSH
Amylase	Cholesterol (Total)	IgA	Phenytoin	Urea
Amylase (Pancreatic)	Cholinesterase	IgG	Phospholipids	Urea Nitrogen
Anti-CCP*	Complement C3	IgM	Phosphorus	Uric Acid
Apolipoprotein A-1	Complement C4	Iron	Potassium	Valproic Acid
Apolipoprotein B	Copper	Iron (TIBC)	Prealbumin	Vancomycin*
ASO*	Cortisol	Iron (UIBC)	Protein Electrophoresis	Vitamin B <sub>12</sub>
AST/SGOT	Creatine Kinase (CK)	Kappa Light Chain*	Protein (Total)	Zinc
Bile Acids*	Creatinine	Lactate Dehydrogenase (LDH)	Quinidine*	
Bilirubin (Direct)	CRP*			

Refer to the package insert of currently available lots for specific analyte and stability claims.  
\*New analytes included in Liquid Assayed Multiquel® Premium.

### Ordering Information

Cat #	Liquid Assayed Multiquel® Premium	Quantity	Cat #	Quantity	
284	Level 1	6 x 5 mL	286	Level 3	6 x 5 mL
285	Level 2	6 x 5 mL	285X	Trilevel MiniPak (1 of each level)	3 x 5 mL



This control is featured in the Unity™ Interlaboratory Program. Participation in the Unity™ Interlaboratory Program can help improve the quality of your laboratory test results and help you meet regulatory and accreditation requirements worldwide. You may also subscribe to one of our Unity™ QC Data Management Solutions designed to improve the effectiveness of your statistical process control. Ask your Bio-Rad Account Representative for more information.

**Discover the power of Unity™ Solutions at [www.QCNet.com](http://www.QCNet.com).**



**Bio-Rad  
Laboratories**

*For further information, please contact the Bio-Rad office nearest you or visit our website at [www.bio-rad.com/chemserumcontrols](http://www.bio-rad.com/chemserumcontrols)*

**Clinical  
Diagnostics Group**

**Website** [www.bio-rad.com/qualitycontrol](http://www.bio-rad.com/qualitycontrol) **Australia** 61-2-9914-2800 **Austria** 43-1-877-8901 **Belgium** 32-9-385-5511 **Brazil** 5521-3237-9400 **Canada** 1-514-334-4372 **China** 86-21-61698500 **Czech Republic** 420-241-430-532 **Denmark** +45-4452-1000 **Finland** 358-9-804-22-00 **France** 33-1-47-95-60-00 **Germany** +49 (0)89-318-840 **Greece** 30-210-7774396 **Hong Kong** 852-2789-3300 **Hungary** +36-1-459-6100 **India** 1800-180-1224 **Israel** 972-3-9636050 **Italy** +39-02-216091 **Japan** 81-3-6361-7070 **Korea** 82-2-3473-4460 **Mexico** +52 (65)5488-7670 **The Netherlands** +31-318-540666 **New Zealand** 64-9-415-2280 **Norway** 47-23-38-41-30 **Poland** 48-22-3319999 **Portugal** 351-21-472-7700 **Russia** 7-495-721-14-04 **Singapore** 65-6415-3170 **South Africa** 27-11-442-85-08 **Spain** 34-91-590-5200 **Sweden** 46-8-555-127-00 **Switzerland** +41 (0)26-674-52-01/02/03 **Taiwan** 886-2-2578-7189 **Thailand** 662-651-8311 **United Kingdom** +44 (0)20-8328-2000